

ZytoLight® SPEC PDGFRA/FIP1L1 TriCheck Probe

For the detection of PDGFRA-FIP1L1 rearrangements by fluorescence *in situ* hybridization (FISH)

The ZytoLight® SPEC PDGFRA/FIP1L1 TriCheck™ Probe is designed to detect rearrangements involving the PDGFRA gene. PDGFRA gene rearrangements most frequently occur in chronic eosinophilic leukemia (CEL), but can be also detected in AML, and T-ALL. The most common gene fusion is FIP1L1-PDGFRA caused by an 800 kb interstitial deletion on chromosome 4q12. Five other partner genes have been identified, including BCR, ETV6, KIF5B, STRN, and CDK5RAP2. Identification of patients harboring a PDGFRA rearrangement is important as these patients respond to a targeted therapy with imatinib. Hence, detection of PDGFRA rearrangements by FISH may be of diagnostic and predictive relevance.

Ordering Information

[Order Online »](#)

ZTV-Z-2209-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Platelet-derived growth factor receptor α

Application FISH

Quantity Sufficient for 5 tests, using 10 μ l probe solution per test.

Technical Info / Product Notes For the Manual please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com